

PRODUCT INFORMATION



Cathelicidin-2 (128-153) (chicken) (trifluoroacetate salt)

Item No. 36947

Formal Name: L-arginyl-L-phenylalanyl-glycyl-L-arginyl-L-phenylalanyl-L-leucyl-L-lysyl-L-isoleucyl-L-arginyl-L-arginyl-L-phenylalanyl-L-arginyl-L-prolyl-L-lysyl-L-valyl-L-threonyl-L-isoleucyl-L-threonyl-L-isoleucyl-L-glutaminyglycyl-L-seryl-L-alanyl-L-arginyl-L-phenylalaninamide, trifluoroacetate salt

Synonyms: CATH-2 (128-153), Fowlicidin-2 (128-153), Myeloid Antimicrobial Peptide 27 (128-153)

Peptide Sequence: RFGRFRLRKIRRFPRPKVTITIQGSARF-NH₂

MF: C₁₄₇H₂₄₅N₅₁O₃₀ • XCF₃COOH

FW: 3,206.9

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

H—Arg—Phe—Gly—Arg—Phe—Leu—Arg—Lys—Ile—Arg—Arg—Phe—Arg—Pro—Lys—Val—Thr—Ile—Thr—Ile—Gln—Gly—Ser—Ala—Arg—Phe—NH₂
• XCF₃COOH

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Cathelicidin-2 (CATH-2) (128-153) (chicken) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the CATH-2 (128-153) in water. We do not recommend storing the aqueous solution for more than one day.

Description

CATH-2 (128-153) is a synthetic antimicrobial peptide that corresponds to amino acids 128-153 of chicken CATH-2.¹ It is active against *E. coli*, *S. aureus*, *S. enteritidis*, and *B. globigii* in a concentration-dependent manner. CATH-2 (128-153) (40 μM) induces hemolysis of isolated chicken erythrocytes but is not cytotoxic to isolated human peripheral blood mononuclear cells (PBMCs). It induces production of chemokine (C-C-motif) ligand 2 (CCL2) and inhibits LPS-induced production of TNF-α, IL-6, IL-8, and IL-10 in isolated human PBMCs.

Reference

- van Dijk, A., Molhoek, E.M., Veldhuizen, E.J.A., et al. Identification of chicken cathelicidin-2 core elements involved in antibacterial and immunomodulatory activities. *Mol. Immunol.* **46(13)**, 2465-2473 (2009).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/10/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM